

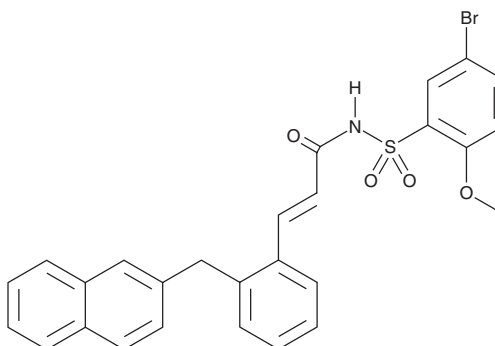
PRODUCT INFORMATION



L-798,106

Item No. 11129

CAS Registry No.: 244101-02-8
Formal Name: (2E)-N-[(5-bromo-2-methoxyphenyl)sulfonyl]-3-[2-(2-naphthalenylmethyl)phenyl]-2-propenamide
Synonyms: CM 9, GW 671021
MF: C₂₇H₂₂BrNO₄S
FW: 536.4
Purity: ≥95%
UV/Vis.: λ_{max}: 221, 285 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

L-798,106 is supplied as a crystalline solid. A stock solution may be made by dissolving the L-798,106 in the solvent of choice, which should be purged with an inert gas. L-798,106 is soluble in organic solvents such as acetonitrile, DMSO, and dimethyl formamide. The solubility of L-798,106 in these solvents is approximately 1, 10, and 20 mg/ml, respectively.

Description

Prostaglandin E₂ (Item No. 14010) exerts its effects through four separate G protein-coupled receptors (EP₁, EP₂, EP₃, and EP₄).¹ L-798,106 is a highly selective EP₃ receptor antagonist with K_i values of 0.3, 916, >5,000, and >5,000 nM at EP₃, EP₄, EP₁, and EP₂, respectively.² At 0.2 μM, it blocks the EP₃ agonist activity of sulprostone (Item No. 14765) on guinea pig vas deferens and trachea.³

References

1. Coleman, R.A., Smith, W.L., and Narumiya, S. Classification of prostanoid receptors: Properties, distribution, and structure of the receptors and their subtypes. *Pharmacol. Rev.* **46**, 205-229 (1994).
2. Juteau, H., Gareau, Y., Labelle, M., et al. Structure-activity relationship of cinnamic acylsulfonamide analogues on the human EP₃ prostanoid receptor. *Bioorg. Med. Chem.* **9**, 1977-1984 (2001).
3. Jones, R.L., Giembycz, M.A., and Woodward, D.F. Prostanoid receptor antagonists: Development strategies and therapeutic applications. *Br. J. Pharmacol.* **158**(1), 104-145 (2009).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/12/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM